

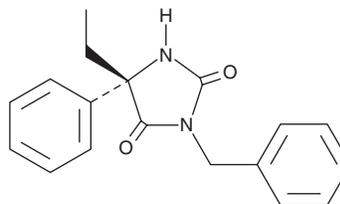
# PRODUCT INFORMATION



## (+)-N-3-Benzyl Nirvanol

Item No. 30097

**CAS Registry No.:** 790676-40-3  
**Formal Name:** (5S)-5-ethyl-5-phenyl-3-(phenylmethyl)-2,4-imidazolidinedione  
**Synonym:** (S)-N-3-benzyl Nirvanol  
**MF:** C<sub>18</sub>H<sub>18</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 294.4  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

(+)-N-3-Benzyl Nirvanol is supplied as a solid. A stock solution may be made by dissolving the (+)-N-3-benzyl Nirvanol in the solvent of choice, which should be purged with an inert gas. (+)-N-3-Benzyl Nirvanol is slightly soluble in DMSO and methanol.

### Description

(+)-N-3-Benzyl Nirvanol is an inhibitor of the cytochrome P450 (CYP) isoform CYP2C19 ( $K_i = 0.25 \mu\text{M}$ ).<sup>1</sup> It is selective for CYP2C19 over CYP1A2, -2A6, -2C8, -2C9, -2D6, -2E1, and -3A4 at 1  $\mu\text{M}$  and over CYP2B6 ( $\text{IC}_{50} = 58 \mu\text{M}$ ).<sup>1,2</sup>

### References

1. Suzuki, H., Kneller, M.B., Haining, R.L., *et al.* (+)-N-3-Benzyl-nirvanol and (-)-N-3-benzyl-phenobarbital: New potent and selective *in vitro* inhibitors of CYP2C19. *Drug Metab. Dispos.* **30(3)**, 235-239 (2002).
2. Walsky, R.L. and Obach, R.S. Verification of the selectivity of (+)N-3-benzyl Nirvanol as a CYP2C19 inhibitor. *Drug Metab. Dispos.* **31(3)**, 343 (2003).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/31/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM